

TECHNICAL SPECIFICATION

CytoFLEX Platform

96-well Plate Sample Loaders



	STANDARD AND DEEP WELL TUBE/PLATE SWITCH CONTROL	STANDARD WELL TUBE/PLATE SWITCH CONTROL	STANDARD WELL MANUAL TUBE/PLATE SWITCH
Part Number, Description	C16574, CytoFLEX Plate Loader DW	C02396, CytoFLEX Plate Loader with Mode Control Upgrade Kit	B63215, CytoFLEX Plate Loader
Tube or Plate Sample Injection Mode	Sample Injection Mode can be changed by using the Sample Injection Mode Control Switch.	Sample Injection Mode can be changed by using the Sample Injection Mode Control Switch.	Sample Injection Mode can be changed manually.
Software	CytExpert version 2.2 or above	CytExpert version 2.0 or above	CytExpert version 1.1 or above
Carryover	<0.5%		
Mixing	Single well mixing. Mixing time is adjustable.		
Plate Compatibility	96-well flat, U-bottom, V-bottom and deep well (45.5 mm height max, 5.8 mm well bottom diameter minimum)	96-well flat, U-bottom and V-bottom	
Minimum Sample Volume	50 µL/well 200 µL/well for efficient deep well mixing	45 µL/well	
Dead Volume	20 µL	20 µL (flat bottom) 10 µL (U- or V-bottom)	
Weight	4.5 kg		
Dimension (W x D x H)	23.5 cm x 25.5 cm x 18.0 cm	23.5 cm x 25.5 cm x 15.5 cm	
Space Requirement	Installed inside the instrument. No extra space is required.		
Power	Connect to the power module of the cytometer directly. No extra power supply is needed.		
Operating Temperature	15-30 °C, non-condensing		
Noise	<65 dBA		

CYTOFLEX AND CYTOFLEX S MODELS

Throughput Standard Well Plates	10 sec acquisition without mixing or backflush: <36 min 10 sec acquisition with 5 sec mixing and 4 sec backflush: <54 min	10 sec acquisition without mixing or backflush: <35 min 10 sec acquisition with 3 sec mixing and 4 sec backflush: <50 min	10 sec acquisition without mixing or backflush: <32 min 10 sec acquisition with 3 sec mixing and 3 sec backflush: <45 min
Throughput Deep Well Plates	10 sec acquisition without mixing and washing: <54 min 10 sec acquisition with 10 sec mixing* and 4 sec backflush: <64 min		

CYTOFLEX LX MODELS

Throughput Standard Well Plates	10 sec acquisition without mixing or backflush: <39 min 5 sec acquisition with 5 sec mixing and 6 sec backflush: <60 min	10 sec acquisition without mixing or backflush: <38 min 10 sec acquisition with 3 sec mixing and 6 sec backflush: <56 min	10 sec acquisition without mixing or backflush: <34 min 10 sec acquisition with 3 sec mixing and 3 sec backflush: <47 min
Throughput Deep Well Plates	10 sec acquisition without mixing and backflush: <40 min 10 sec acquisition with 6 sec mixing and 4 sec backflush: <69 min		

*Performance with deep well plates are based on the testing results with Beckman Coulter deep well plates, part number 267001, 267006 and 140504. Results may vary with different types of plates.



For research use only. Not for use in diagnostic procedures.
© 2020 Beckman Coulter Life Sciences. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at beckman.com

FLOW-3308SPEC04.20